

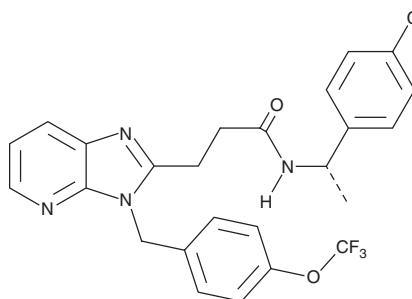
PRODUCT INFORMATION



CRT0273750

Item No. 20904

CAS Registry No.: 1979939-16-6
Formal Name: N-[(1S)-1-(4-chlorophenyl)ethyl]-3-[[4-(trifluoromethoxy)phenyl]methyl]-3H-imidazo[4,5-b]pyridine-2-propanamide
MF: C₂₅H₂₂ClF₃N₄O₂
FW: 502.9
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

CRT0273750 is supplied as a solid. A stock solution may be made by dissolving the CRT0273750 in the solvent of choice, which should be purged with an inert gas. CRT0273750 is soluble in DMSO.

Description

CRT0273750 is an ectonucleotide pyrophosphatase/phosphodiesterase family member 2 (ENPP2) inhibitor (IC₅₀ = 1 nM).¹ It prevents lysophosphatidylcholine hydrolysis in a plasma choline release assay (IC₅₀ = 13 nM). CRT0273750 inhibits the migration of 4T1 cells in a migration assay (IC₅₀ = 24 nM). It decreases plasma levels of the ENPP2 hydrolysis product lysophosphatidic acid in an MDA-MB-231 mouse xenograft model.

Reference

1. Shah, P., A., C., Foxton, C., *et al.* Discovery of potent inhibitors of the lysophospholipase autotaxin. *Bioorg. Med. Chem. Lett.* **26(22)**, 5403-5410 (2016).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/16/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM